

---

## DGK $\eta$ (DGKH)

Product code 12-406-20N

Catalytic domain of human DGK $\eta$  [1-1147 amino acids of accession number NP\_821077.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (131 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform : ADP-Glo  
Substrate : Diacylglycerol

ATP ( $\mu$ M) Km app / Bin : 24 / 25  
Metal : Mg  
Reference compound : Non-disclosable  
IC50 at ATP Bin (nM) :  
IC50 at 1 mM ATP (nM) :

---